

Material identification

Name: FlexMAP™ Microspheres
Product #: L100-Uxxx-01 and 14-00xxx

Composition

Hazards identified with this product are those associated with the following components.

Ingredient:	Concentration (% wt):
Proprietary dyes	< 1%
Anti-tag Oligonucleotides	< 0.01%
xMAP Carboxylated Microspheres	< 1%
Tris- EDTA buffer (10mM Tris, 1mM EDTA)	> 98%

Hazards identification

- Irritating to eyes, respiratory system, and skin.
- May be toxic in contact with skin and if swallowed.
- In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

First aid measures

- Wear suitable protective clothing.
- If eye contact, flush eyes with copious amounts of water for 15 minutes. Assure adequate flushing by separating the eyelids with fingers. Call a physician.
- If skin contact, immediately wash skin with soap and copious amounts of water.
- If swallowed, wash out mouth with water, provided person is conscious. Call a physician.
- If inhaled, remove to fresh air. If not breathing, give artificial respiration. If breathing is difficult, give oxygen.
- Wash contaminated clothing before reuse.

Fire and explosion hazard data

Extinguishing media: Water, foam, dry chemical, or CO₂.

Special fire-fighting procedures: Wear self-contained breathing apparatus and protective clothing to prevent contact with skin and eyes.

Unusual fire and explosion hazards: Emits toxic fumes under fire conditions.

Accidental release measures/disposal considerations

Steps to be taken in case material is released or spilled:

- Wear respirator, chemical safety goggles or face shield, and other protective clothing such as rubber boots and heavy rubber gloves.
- Take up with absorbent material. Avoid raising dust. Place in a bag and hold for waste disposal.
- Ventilate area and wash spill site after material pick-up is complete.

Waste disposal method: Comply with all applicable federal, state, or local regulations.

Handling and storage

Store in sealed container in cool (2-8°C), dry location. Keep away from oxidizers. See Exposure controls and personal protection section below.

Exposure controls and personal protection

- Wear appropriate respirator, chemical-resistant gloves, safety goggles or face shield, and other protective clothing.
- Safety shower and eye bath.
- Irritant.
- Avoid contact with eyes, skin, and clothing.
- Wash thoroughly after handling.
- Keep tightly closed.

Stability and reactivity

Stability: Stable.

Incompatibilities: Strong oxidizing agents.

Hazardous decomposition products: Toxic fumes of carbon monoxide, carbon dioxide, nitrogen oxides and sulfur oxides.

Hazardous polymerization: Has not been reported.

Toxicological information

Acute effects: May be harmful by inhalation, ingestion, or skin absorption; causes eye and skin irritation, material is irritating to mucus membranes and upper respiratory tract.

To our knowledge, the chemical, physical, and toxicological properties have not been thoroughly investigated.

Ecological information

Data not yet available.

Other information

The above information is believed to be correct but is not all-inclusive and should only be used as a guide. Luminex Corporation will not be held liable for any damage resulting from handling or contact with the above product if used, handled, stored, or disposed of in a manner other than as prescribed.